

JARID1A Protein

(Catalog No. E24020)

Description

Human JARID1A, also known as KDM5A and RBBP2, GenBank Accession #NM_001042603), a.a. 1-1090 with C-terminal FLAG-tag

Molecular Weight (kDa)

125 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 80 µg/ml FLAG peptide, and 20% glycerol

Source

Expressed in Sf9 cells via a baculovirus expression system

Size

20 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity > 52%

Application Notes

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Ordering Information:

Products	Size	Cat. No.
JARID1A Protein	20 µg	E24020 -1

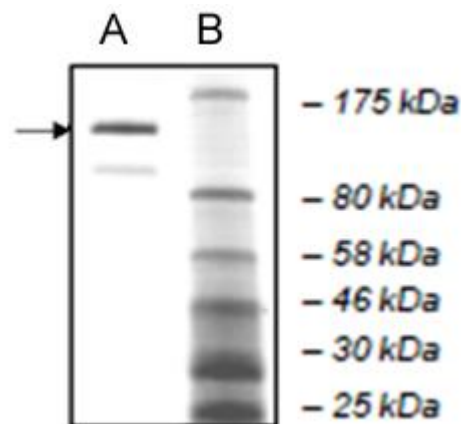


Fig. 1. 4-20% SDS-page Coomassie staining.
A: JARID1A 2 µg; B: Marker

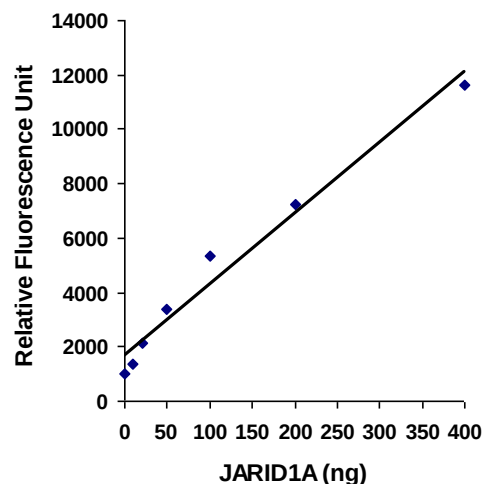


Fig. 2. Activity of recombinant JARID1A protein was measured with the **Epigenase JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric)** (Cat. #P-3083)

This product is for research purposes only. Not intended for use in diagnostic procedures.